

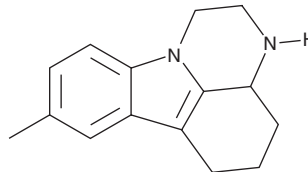
PRODUCT INFORMATION



Pirlindole

Item No. 22060

CAS Registry No.: 60762-57-4
Formal Name: 2,3,3a,4,5,6-hexahydro-8-methyl-1H-pyrazino[3,2,1-jk]carbazole
MF: C₁₅H₁₈N₂
FW: 226.3
Purity: ≥98%
UV/Vis.: λ_{max}: 229, 275 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Pirlindole is supplied as a crystalline solid. A stock solution may be made by dissolving the pirlindole in the solvent of choice. Pirlindole is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of pirlindole in these solvents is approximately 5 and 1 mg/ml, respectively.

Description

Pirlindole is a selective and reversible monoamine oxidase A (MAO-A) inhibitor (IC₅₀s = 250 and 34.2 nM for rat brain and heart MAO-A, respectively).¹ It is selective for MAO-A over MAO-B (K_is = 52,100 and 59,900 nM, for rat brain and heart MAO-B, respectively). In rats, it reverses the depressive-like effects induced by chronic mild stress (CMS), increases proliferation of hippocampal neural progenitor cells, and reverses dendritic atrophy in granule neurons.² Pirlindole is also an inhibitor of enterovirus-D68 and coxsackievirus B3 (CV-B3), inhibiting the genome replication phase of CV-B3 infection with an EC₅₀ value of 7.7 μM independent of MAO-A activity.³

References

1. Bruhwyler, J., Liégeois, J.F., and Géczy, J. Pirlindole: A selective reversible inhibitor of monoamine oxidase A. A review of its preclinical properties. *Pharmacol. Res.* **36(1)**, 23-33 (1997).
2. Morais, M., Santos, P.A., Mateus-Pinheiro, A., *et al.* The effects of chronic stress on hippocampal adult neurogenesis and dendritic plasticity are reversed by selective MAO-A inhibition. *J. Psychopharmacol.* **28(12)**, 1178-1183 (2014).
3. Ulferts, R., de Boer, S.M., van der Linden, L., *et al.* Screening of a library of FDA-approved drugs identifies several enterovirus replication inhibitors that target viral protein 2C. *Antimicrob. Agents Chemother.* **60(5)**, 2627-2638 (2016).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/22/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM